

## **Product Data Sheet**

Product Name: Scrambled-β-Amyloid (1-42), 5-FAM-labeled

Catalog Number: 60892 (0.1 mg) Lot Number: See label on vial

Sequence: 5-FAM-Ala-Ile-Ala-Glu-Gly-Asp-Ser-His-Val-Leu-Lys-Glu-Gly-Ala-Tyr-

Met-Glu-Ile-Phe-Asp-Val-Gln-Gly-His-Val-Phe-Gly-Gly-Lys-Ile-Phe-Arg-

Val-Val-Asp-Leu-Gly-Ser-His-Asn-Val-Ala-OH (3-letter code)

5-FAM-AIAEGDSHVLKEGAYMEIFDVQGHVFGGKIFRVVDLGSHNVA

(1-letter code)

Molecular Weight: 4873.4

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Using basic buffer (1%NH<sub>4</sub>OH), reconstitute by adding 50 μl to 0.1 mg Scrambled-β-Amyloid (1-42), 5-FAM-labeled peptide. Dilute this stock solution with working buffer, such as PBS or other buffers.

Storage: Scrambled- $\beta$ -Amyloid (1-42), 5-FAM-labeled peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at  $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}$ C or lower.

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Scrambled A $\beta$ 42 (Anaspec, San Jose, CA, USA) was also spotted to determine background, nonspecific binding values for each probe. Radioactivity was quantified in a scintillation counter. The amount of specific binding was calculated by subtracting the radioactivity of the scrambled A $\beta$ 42 spot from the A $\beta$ 42 spot for each probe. Shaked, BM. The FASEB Journal. 208, 1254 (2006).

## **Published Citations:**

Shaked, BM. The FASEB Journal. 208, 1254 (2006).

## **Related Products:**

NameCat #Sizeβ-Amyloid (1-42), FAM-labeled23526-010.1 mgFAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA23525-050.5 mg

β-Amyloid (1-42), TAMRA-labeled TAMRA-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	60476	0.1 mg
β-Amyloid (1-42), AMCA-LC-labeled AMCA-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	60475-01	0.1 mg
β-Amyloid (1-42), Sulforhodamine 101-labeled Sulforhodamine 101-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMV	62221 GGVVIA	0.1 mg
β-Amyloid (1-42), HiLyte Fluor <sup>TM</sup> 488 - labeled HiLyte Fluor <sup>TM</sup> 488-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVG	60479-01 GVVIA	0.1 mg
β-Amyloid (1-42), HiLyte Fluor™ 555-labeled HiLyte Fluor 555-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGG	60480-01 VVIA	0.1 mg

For Research Use Only